



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

HDAC 6

human, recombinant, N-terminal GST tag
Catalog #: 50006

Formulated in: 25 mM Tris-HCl, pH 8.0,
138 mM NaCl, 0.05% Tween-20 and 10%
glycerol.

Stability: >6 months at -80°C

Reference:

1. Strausberg, R.L. et al., Proc. Natl. Acad. Sci.
U.S.A. 99 (26), 16899-16903 (2002).

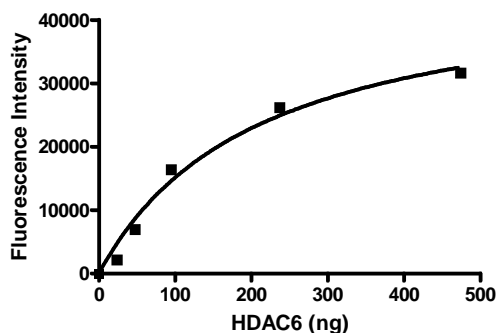
Description: Human HDAC6 (GenBank
Accession No. BC069243), full length with
N-terminal GST tag, MW= 159 kDa,
expressed in baculovirus expression system.

Specific Activity: 257 U/μg. One U =1
pmol/min, Assay condition: 25 mM Tris/Cl,
pH8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM
MgCl₂, and 0.1 mg/ml BSA, 30 μM BPS
HDAC substrate (Catalog number 50037),
and 5 ng/μl HDAC6. Incubation condition:
30 min at 37°C.

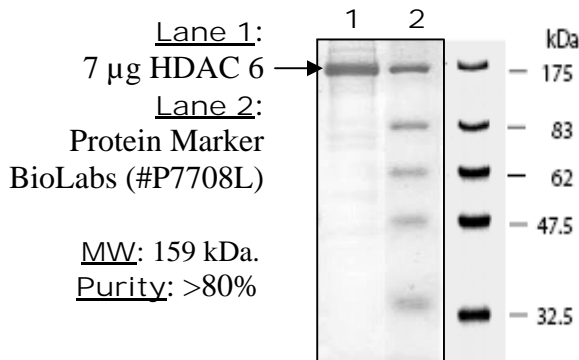
Application: Useful for the study of
enzyme kinetics, screening inhibitors, and
selectivity profiling.

Quality Assurance

HDAC6 Specific Activity



10% SDS-PAGE
Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Assay Protocol

Material: Assay buffer (BPS catalog number 50031); Assay developer (BPS catalog number 50030); HDAC Substrate (BPS number 50032)

Step 1: adding all reaction mixture to a low binding black plate

35 µl of HDAC assay buffer (BPS catalog number 50031)
5 µl of 1 mg/ml BSA
5 µl of 200 uM substrate (BPS catalog number: 50037)
5 µl of HDAC6 (0.01 ug/ul)

Always add HDAC at the last.

Incubate at 37 °C for 30 min.

Setp2: stop the reaction

add 50 µl of fluor de lys developer (2x) (BPS catalog number 50030) and incubate the plate at room temperature for 15 min

Step 3: read sample in a microtiter-plate reading fluorimeter capable of excitation at a wavelength in the range 350-380 nm and detection of emitted light in the range 440-460 nm.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com